

## SFX (6-(Fluorescein-6-carboxamido)hexanoic acid) Succinimidyl Ester, 5 Isomer



### Product attributes

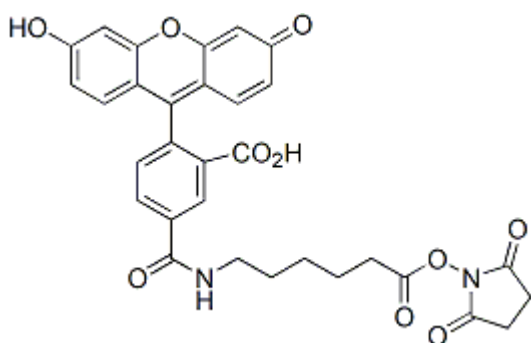
Excitation/Emission	494/521 nm (pH 9)
---------------------	-------------------

5-SFX has a 6-atom aminohexanoyl spacer group that separates the fluorescein dye and the amine-reactive group.

### Product Description

5-SFX (full name: 6-(Fluorescein-5-carboxamido)hexanoic acid, succinimidyl ester, single isomer) has a 6-atom aminohexanoyl spacer group that separates the fluorescein dye and the amine-reactive group.

- $\lambda_{Ex}/\lambda_{Em}$  (pH 9) = 494/521 nm
- Yellow solid soluble in DMF, DMSO or pH>6.5 buffers
- Store at -20 °C and protect from light
- $C_{31}H_{26}N_2O_{10}$
- MW:586.55



This datasheet was generated on June 9, 2026 at 03:34:45 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/5-sfx-6-fluorescein-5-carboxamido-hexanoic-acid-succinimidyl-ester-single-isomer/>